

Compatibility of Single Step (KRX) Competent Cells with the MagneGST™ Pull-Down System

ABSTRACT

This study demonstrates the use of Single Step (KRX) Competent Cells for studying protein:protein interactions using the MagneGST™ Pull-Down System. GST-fusion bait protein expressed in KRX was captured on MagneGST™ glutathione-linked magnetic particles and incubated with prey protein expressed in vitro in the TNT® T7 Quick Coupled Transcription/Translation System. Captured prey was detected using HaloTag® Interchangeable Labeling Technology.

Promega Notes 96, 22–23.

Trista Schagat

Promega Corporation

Publication Date: 2007



Download Article

(127 KB)
